p53 (human), (recombinant) (GSTtag)

p53 is a much studied and complex multifunctional protein, which acts as a tumor suppressor in many tumor types and induces growth arrest or apoptosis depending on the physiological circumstances and cell type. It is involved in cell cycle regulation as a trans-activator that acts to negatively regulate cell division by controlling a set of genes required for this process. One of the activated genes is an inhibitor of cyclin-dependent kinases. Apoptosis induction appears to be mediated either by stimulation of BAX and FAS antigen expression, or by repression of Bcl-2 expression. The p53 pathway responds to stresses that can disrupt the fidelity of DNA replication and cell division. A stress signal is transmitted to the p53 protein by post-translational modifications. Such modification results in the activation of p53 protein as a transcription factor that initiates a program of cell cycle arrest, cellular senescence or apoptosis.

Citations: 4

View Online »

Ordering Information

Order Online »

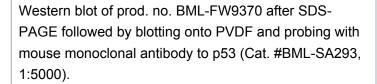
BML-FW9370-0050 5

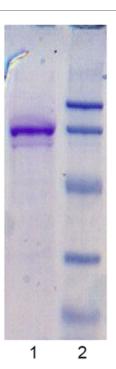
50µg

Manuals, SDS & CofA

View Online »







Coomassie stained gel of prod.no. BML-FW9370: Lane 1: FW9370 (3µg), Lane 2: MW markers (top to bottom) 113, 93, 50, 35, and 28 kDa.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

TRP53, Tumor protein p53 **Alternative Name**

Formulation Liquid. In 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM

DTT, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 25%

glycerol.

MW ~70kDa (~85/90kDa observed with GST tag)

Purity ≥90% (SDS-PAGE)

Source Produced in E. coli. Full length human p53 is fused to a

GST-tag.

UniProt ID P04637

Last modified: September 18, 2025



eu@enzolifesciences.com